

# YC-1

## Guanylyl cyclase activator

Nitric oxide (NO) independent, superoxide-sensitive stimulator of soluble guanylyl cyclase (sGC),  $EC_{50}=18.6\mu M$ . Induces concentration-dependent increase in cGMP levels in cultured cells and in rabbit aortic rings. Inhibits platelet adhesion to collagen. Non-specific phosphodiesterase inhibitor. YC-1 inhibits HIF- $\alpha$  activity in tumors resulting in blocked angiogenesis and an inhibition of tumor growth. Inhibits  $Na^+$  channel. Binds to [ $^3$ H]BW202W92.

Citations: 15

[View Online »](#)

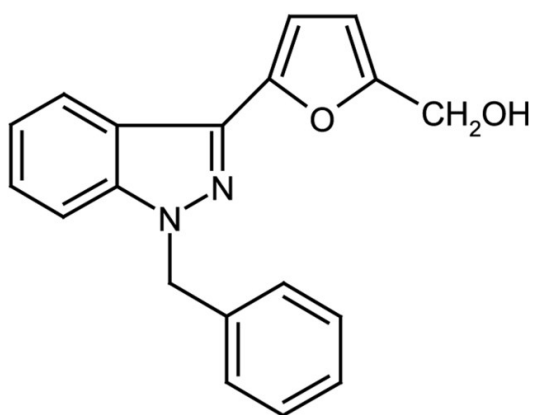
## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-420-025-M001 | 1mg |
| ALX-420-025-M005 | 5mg |

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |                                                                  |
|-------------------|------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. |
| Handling          | Protect from light.                                              |
| Long Term Storage | +4°C                                                             |
| Shipping          | Ambient Temperature                                              |

## Regulatory Status

RUO - Research Use Only

## Product Details

|               |                                               |
|---------------|-----------------------------------------------|
| Appearance    | White to pale yellow powder.                  |
| CAS           | 170632-47-0                                   |
| Couple Target | Guanylyl cyclase                              |
| Couple Type   | Activator                                     |
| Formula       | $C_{19}H_{16}N_2O_2$                          |
| Identity      | Identity determined by 1H-NMR.                |
| MW            | 304.4                                         |
| Purity        | ≥98% (HPLC)                                   |
| Solubility    | Soluble in DMSO, methanol or dichloromethane. |

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)